

DMRIE-C Reagent

Part no. 10459.pps MAN0000829 Rev. Date 14 July 2011

Cat. no. 10459-014 Size 1 mL Store at 4°C (do not freeze)

Description

DMRIE-C Reagent is suitable for transfecting DNA and RNA into eukaryotic cells, and is particularly effective for transfecting suspension cells (e.g. Jurkat) and other lymphoid-derived cell lines. Refer to the Cell Lines database at www.invitrogen.com for a list of cell types successfully transfected. DMRIE-C can also be used for *in vivo* delivery of DNA. DMRIE-C is a 1:1 (M/M) liposome formulation of the cationic lipid DMRIE (1,2-dimyristyloxy-propyl-3-dimethyl-hydroxy ethyl ammonium bromide) and cholesterol in membrane filtered water.

Important Guidelines for Transfection

- Form complexes using the recommended amounts of DNA (or RNA) and DMRIE-C. Optimize as necessary.
Note: We recommend diluting DNA (or RNA) and DMRIE-C into Opti-MEM® I Reduced Serum Medium (Cat. no. 31985-062) before complexing.
- DMRIE-C is a lipid suspension that may settle with time. Mix thoroughly by inverting the tube 5–10 times before use.
- Transfect cells at the recommended confluence or cell density. Optimize as necessary. Maintain the same seeding conditions between experiments.
- *Do not* add antibiotics to media during transfection.
- For optimal results, perform transfection in medium without serum. Cells may be transfected in the presence of serum, if desired. However, complexes *must* be formed in serum-free medium.
- Test serum-free media for compatibility with DMRIE-C since some serum-free formulations (e.g. CD 293, 293 SFM II, VP-SFM) may inhibit cationic lipid-mediated transfection.

Intended Use: For research use only.

Not intended for any animal or human therapeutic or diagnostic use.

Transfecting Adherent Mammalian Cells with DNA

Use the following procedure to transiently or stably transfect DNA into mammalian cells. All amounts and volumes are given on a *per-well basis*.

1. One day before transfection, plate cells in growth medium without antibiotics such that they will be at the recommended confluence at the time of transfection.

Condition	Cell no.	Growth med. vol.	Format	Confluence at time of transfection
Transient	$1-2 \times 10^5$	2 mL	6-well	60–80%
Stable	$1-2 \times 10^5$	4 mL	60-mm	30–50%

2. For each transfection sample, prepare complexes as follows:
 - a. Dilute 1–2 μg of DNA in 500 μL (1 mL for stable transfection) of Opti-MEM[®] I Reduced Serum Medium (or other medium) without serum.
 - b. Mix DMRIE-C before use, then dilute 2–12 μL of DMRIE-C in 500 μL (1 mL for stable transfection) of Opti-MEM[®] I Medium (or other medium) without serum. Let stand at room temperature for 30–45 minutes.
 - c. Combine the diluted DNA with diluted DMRIE-C (total volume = 1 mL; 2 mL for stable transfection). Mix gently and incubate for 15–45 minutes at room temperature (the solution may appear cloudy).
3. Remove the growth medium from the cells and wash once with 2 mL of growth medium without serum. Remove the wash medium.
4. Add the complexes (from step 2c of this procedure) to cells. Mix gently by rocking the plate back and forth.
5. Incubate cells at 37°C in a CO₂ incubator for 4–24 hours.
6. Replace the medium with 2 mL of growth medium containing serum (4 mL for stable transfection). Alternatively, add 1 mL of growth medium containing 2X the normal concentration of serum (2 mL for stable transfection) without removing the DNA-containing medium.
7. **Transient:** Test for transgene expression 24–72 hours post-transfection.
Stable cell lines: Passage cells at a 1:5 (or higher dilution) into selective medium 48–72 hours post-transfection.

Transfecting Suspension Mammalian Cells with DNA

Use the following procedure to transfect mammalian cells in suspension in a *6-well format*. All amounts and volumes are given on a *per-well basis*.

1. *For each transfection sample*, prepare complexes as follows:
 - a. Aliquot 500 μL of Opti-MEM[®] I Reduced Serum Medium (or other medium) without serum into a well. Mix DMRIE-C before use, then dilute 2–12 μL into the medium, and mix gently by swirling the plate.
 - b. In a 12 \times 75 mm polystyrene tube, dilute 4 μg of DNA in 500 μL of Opti-MEM[®] I Medium (or other medium) without serum.
 - c. Add the diluted DNA to the well containing diluted DMRIE-C (total volume = 1 mL). Mix by swirling the plate and incubate for 15–45 minutes at room temperature (solution may appear cloudy).
2. While complexes are forming, prepare a single-cell suspension from stock cells. Wash the cells once with serum-free growth medium without antibiotics, and add $2\text{--}3 \times 10^6$ cells in 0.2 mL of serum-free growth medium without antibiotics to each well.
3. Incubate cells at 37°C in a CO₂ incubator for 4–5 hours.
4. Add 2 mL of growth medium containing 15% fetal bovine serum to the cells. **Note:** For Jurkat and MOLT-4 cells, add 50 $\mu\text{g}/\text{mL}$ PMA and 1 $\mu\text{g}/\text{mL}$ PHA-L, if desired, to enhance promoter activity and gene expression. For K562 and KG-1 cells, add PMA alone.
5. Test for transgene expression 24–48 hours post-transfection.

Transfecting Mammalian Cells with RNA

Use the following procedure to transfect mammalian cells with RNA in a *6-well format*. All amounts and volumes are given on a *per-well basis*.

1. One day before transfection, plate $2\text{--}3 \times 10^5$ cells in 2 mL of growth medium without antibiotics such that they will be 80% confluent at the time of transfection.
2. Just before transfection, remove the growth medium from the cells and wash once with 2 mL of growth medium without serum. Remove wash medium.
3. *For each transfection sample*, prepare complexes as follows:

To a 12 \times 75 mm polystyrene tube, add 1 mL of Opti-MEM[®] I Medium (or other medium) without serum. Mix DMRIE-C, then dilute 2–12 μL into the medium. Mix or vortex briefly. Proceed to step 4 on page 4.

Transfecting Mammalian Cells with RNA, Continued

4. For each transfection sample, add 2–5 μg of RNA into the tube containing DMRIE-C and medium (from step 3 of this procedure, page 3). Mix or vortex briefly, then immediately add complexes to washed cells.
5. Incubate at 37°C in a CO₂ incubator for 4 hours, then replace transfection medium with complete growth medium.
6. Assay for expression of transfected RNA 16–24 hours post-transfection.
Note: Capped and polyadenylated mRNA is translated more efficiently and is more stable within the cell.

Optimizing Transfection

To obtain the highest transfection efficiency and low non-specific effects, optimize transfection conditions by varying cell density, DNA (or RNA) and DMRIE-C concentrations, and transfection incubation time.

Scaling Up or Down Transfections

Procedures are provided to transfect cells in a 6-well format (60-mm format for stable transfection with DNA). For other formats, vary the amounts of DNA (or RNA), DMRIE-C Reagent, cells, and medium used in proportion to the relative surface area of the tissue culture vessel.

Certificate of Analysis

The Certificate of Analysis provides detailed quality control information for each product. Certificates of Analysis are available on our website. Go to www.lifetechnologies.com/support and search for the Certificate of Analysis by product lot number, which is printed on the box.

Limited Use Label License

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

©2011 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

For support visit www.lifetechnologies.com/support or email techsupport@lifetech.com

www.lifetechnologies.com

The logo for Life Technologies, featuring the word "life" in a lowercase, cursive script font, with "technologies™" in a smaller, sans-serif font below it.